

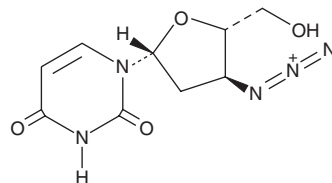
PRODUCT INFORMATION



3'-Azido-2',3'-dideoxyuridine

Item No. 36544

CAS Registry No.: 84472-85-5
Formal Name: 3'-azido-2',3'-dideoxy-uridine
Synonyms: AZddU, AZU, Navuridine, NSC 380882
MF: $C_9H_{11}N_5O_4$
FW: 253.2
Purity: $\geq 95\%$
UV/Vis.: $\lambda_{\max}$: 261 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

3'-Azido-2',3'-dideoxyuridine (AZddU) is supplied as a solid. A stock solution may be made by dissolving the AZddU in the solvent of choice, which should be purged with an inert gas. AZddU is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AZddU in these solvents is approximately 2 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of AZddU can be prepared by directly dissolving the solid in aqueous buffers. The solubility of AZddU in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

AZddU is an antiviral nucleoside analog.¹ It is phosphorylated by thymidine kinase.² AZddU inhibits cytopathogenicity induced by HIV or Moloney sarcoma virus (MSV) in MT-4 cells (EC_{50} s = 0.36 and 15 μM , respectively).¹

References

1. Balzarini, J., Baba, M., Pauwels, R., *et al.* Anti-retrovirus activity of 3'-fluoro- and 3'-azido-substituted pyrimidine 2',3'-dideoxynucleoside analogues. *Biochem. Pharmacol.* **37(14)**, 2847-2856 (1988).
2. Eriksson, B.F., Chu, C.K., and Schinazi, R.F. Phosphorylation of 3'-azido-2',3'-dideoxyuridine and preferential inhibition of human and simian immunodeficiency virus reverse transcriptases by its 5'-triphosphate. *Antimicrob. Agents Chemother.* **33(10)**, 1729-1734 (1989).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/01/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM